

PRODUCT INFORMATION



MK-571 (sodium salt)

Item No. 70720

CAS Registry No.: 115103-85-0

Formal Name: (E)-3-[[[3-[2-(7-chloro-2-quinolinyl)ethenyl]phenyl][3-(dimethylamino)-3-oxopropyl]thio]methyl]thio]-propanoic acid, sodium salt

Synonym: L-660,711

MF: $C_{26}H_{26}ClN_2O_3S_2 \cdot Na$

FW: 537.1

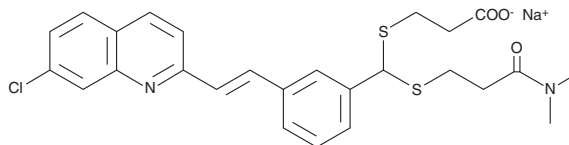
Purity: $\geq 95\%$

UV/Vis.: λ_{max} : 226, 283, 328, 345, 358 nm

Supplied as: A crystalline solid

Storage: $-20^\circ C$

Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

MK-571 (sodium salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the MK-571 (sodium salt) in the solvent of choice, which should be purged with an inert gas. MK-571 (sodium salt) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). It is also soluble in water. The solubility of MK-571 (sodium salt) in ethanol is approximately 1 mg/ml, approximately 10 mg/ml in DMSO and DMF, and approximately 20 mg/ml in water. We do not recommend storing the aqueous solution for more than one day.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of MK-571 (sodium salt) can be prepared by directly dissolving the crystalline compound in aqueous buffers. The solubility of MK-571 (sodium salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

MK-571 is a cysteinyl leukotriene 1 (CysLT₁) receptor antagonist ($K_i = 2.1$ nM in a radioligand binding assay using isolated human lung membranes).¹ It inhibits contractions induced by leukotriene D₄ (LTD₄; Item No. 20310) or LTE₄ (Item No. 20410) in histamine-primed isolated guinea pig trachea ($pA_{2S} = 9.4$ and 9.1 , respectively), but does not inhibit contractions induced by LTC₄ (Item No. 20210) in histamine-primed isolated guinea pig trachea when used at a concentration of 190 nM. MK-571 (100 nM) inhibits LTD₄-induced calcium mobilization in COS-7 monkey kidney cells expressing the human CysLT₁ receptor in a reporter assay.² MK-571 (5 mg/kg) improves tissue damping and elasticity, markers of lung function, and decreases IL-4 and IL-5 levels in bronchoalveolar lavage fluid (BALF) in a mouse model of ovalbumin-induced asthma.³

References

1. Jones, T.R., Zamboni, R., Belley, M., et al. *Can. J. Physiol. Pharmacol.* **67**(1), 17-28 (1989).
2. Lynch, K.R., O'Neill, G.P., Liu, Q., et al. *Nature* **399**(6738), 789-793 (1999).
3. da Cunha, A.A., Silveira, J.S., Antunes, G.L., et al. *Exp. Lung Res.* **47**(8), 355-367 (2021).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/06/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM